

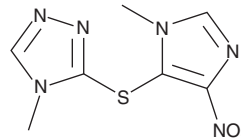
PRODUCT INFORMATION



MNITMT

Item No. 43554

CAS Registry No.: 177653-76-8
Formal Name: 4-methyl-3-[(1-methyl-4-nitro-1H-imidazol-5-yl)thio]-4H-1,2,4-triazole
Synonym: NSC 631156
MF: C₇H₈N₆O₂S
FW: 240.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MNITMT is supplied as a solid. A stock solution may be made by dissolving the MNITMT in the solvent of choice, which should be purged with an inert gas. MNITMT is slightly soluble (0.1-1 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

MNITMT is an immunosuppressant.¹ It inhibits cell proliferation in the human mixed lymphocyte reaction (EC₅₀ = 2.7 μM). MNITMT (45 mg/kg per day) increases allograft survival in a mouse skin transplantation model. It also inhibits the human IgG-induced antibody response in rabbits when administered at a dose of 2 mg/kg per day.²

References

1. Crawford, D.J., Maddocks, J.L., Jones, D.N., *et al.* Rational design of novel immunosuppressive drugs: Analogues of azathioprine lacking the 6-mercaptopurine substituent retain or have enhanced immunosuppressive effects. *J. Med. Chem.* **39**(14), 2690-2695 (1996).
2. Al-Safi, S., Tashtoush, B., and Hassan, M. Comparison of the effects of azathioprine and its novel non-mercaptopurine analog on antibody response in rabbits. *Pol. J. Pharmacol.* **55**(2), 239-243 (2003).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/04/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM